

Blocking peptide for Adiponectin receptor 1 pAb (Prod. No. ALX- 210-645)

Synthetic peptide corresponding to aa 345-369 of human, mouse or rat adiponectin receptor 1 (AdipoR1).

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-151-045-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Some of the peptide may precipitate upon thawing. Resuspend by sonification and dilute with appropriate buffer to dissolve if necessary.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Concentration	3.5mg/ml
Formulation	Liquid. In 20mM acetate buffer, pH 4.5.
Purity	≥90% (HPLC)
Sequence	V ³⁴⁵ AAAFVHFYGVSNLQEFRYGLEGGC ³⁶⁹

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com